

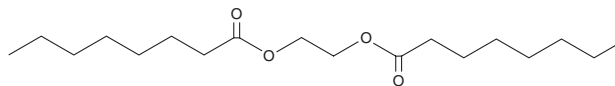
PRODUCT INFORMATION



Diocanoyl Glycol

Item No. 31515

CAS Registry No.: 627-86-1
Formal Name: octanoic acid,
1,1'-(1,2-ethanediyl) ester
Synonyms: Diocanoyl Glycol,
diC8-Ethylene Glycol,
Ethylene Glycol Diocanoate,
8:0 Ethylene Glycol, NSC 97300,
SFE 18:1;O2



MF: C₁₈H₃₄O₄
FW: 314.5
Purity: ≥98%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Diocanoyl glycol is supplied as a neat oil. A stock solution may be made by dissolving the diocanoyl glycol in the solvent of choice, which should be purged with an inert gas. Diocanoyl glycol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of diocanoyl glycol in these solvents is approximately 30 mg/ml.

Description

Diocanoyl glycol is a diacylglycerol kinase (DGK) inhibitor ($K_i = 58 \mu\text{M}$).¹ It reduces basal and thrombin-stimulated phosphatidic acid levels in washed isolated human platelets when used at concentrations of 30, 100, and 500 μM .

References

1. Bishop, W.R., Ganong, B.R., and Bell, R.M. Attenuation of sn-1,2-diacylglycerol second messengers by diacylglycerol kinase. Inhibition by diacylglycerol analogs in vitro and in human platelets. *J. Biol. Chem.* **261**(15), 6993-7000 (1986).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/14/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM